

Anti-PYK2

Rabbit Polyclonal Antibody

Catalog # P92-63R

Lot # M132-7

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the PYK2 (FAK2) protein

Cross Reactivity

- Western blot human PYK2

PYK2 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 400-404 of human PYK2

Formulation

PBS, pH7.4, 0.02%NaN₃, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

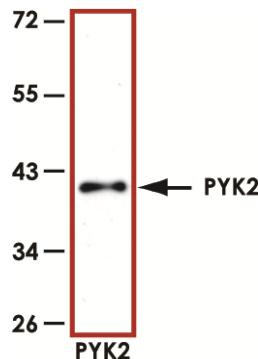
Scientific Background

PYK2 (also known as FAK2/RAFTK) is a member of the focal adhesion PTK family. PYK2/FAK2 can be activated by a variety of extracellular signals that elevate intracellular calcium concentration, and by stress signals (1). Unlike FAK, which is widely expressed in various tissues and links transmembrane integrin receptors to intracellular pathways, PYK2/FAK2 is expressed mainly in the central nervous system and in cells derived from hematopoietic lineages. In osteoclasts, although FAK is expressed, PYK2/FAK2 appears to be the predominant mediator of integrin alpha(v)beta3 signaling events that influence osteoclast physiology and pathology (2).

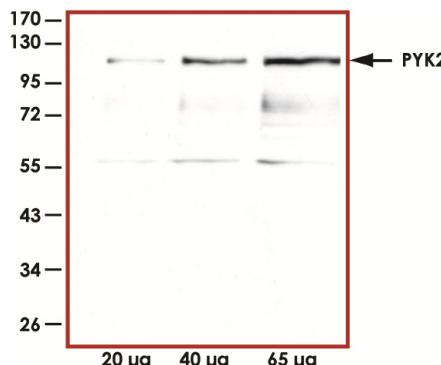
References

1. Avraham, H. et al: RAFTK/Pyk2-mediated cellular signalling. *Cell Signal.* 2000 Mar;12(3):123-33.
2. Xiong, W C. et al: PYK2 and FAK in osteoclasts. *Front Biosci.* 2003 Sep 1;8:d1219-26.

Sample Data



Representative Western blot with Anti-PYK2 (1:1000) using 2 ng of His-tagged human recombinant PYK2.



Representative Western blot with Anti-PYK2 (1:1000) using 20 µg, 40 µg and 65 µg of Jurkat cell lysate.

Anti-PYK2

Rabbit Polyclonal Antibody

Catalog Number	P92-63R
Specific Lot Number	M132-7
Purification	Affinity chromatography
Concentration	1.0 µg/µL
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.